

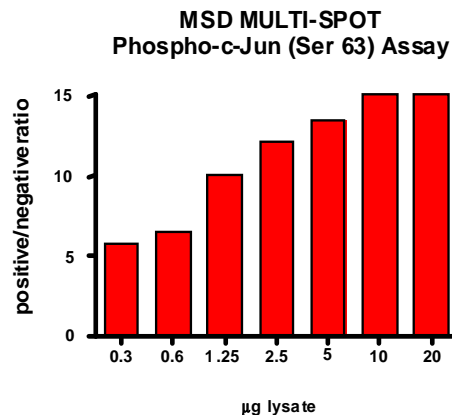
Meso Scale Discovery® Whole Cell Lysate Set

Phospho-c-Jun (Ser 63)

Catalog No:	C11CG-1
Contents:	2 x 100 µg MSDLY0026 pc-Jun Negative Cell Lysate Cell lysate from serum deprived NIH3T3 cells 2 x 100 µg MSDLY0061 pc-Jun Positive Cell Lysate Cell lysate from NIH3T3 cells harvested 15 minutes after UV irradiation (40 mJ/cm ²) to stimulate c-Jun phosphorylation
Concentration:	2 mg/mL in MSD® Complete Tris Lysis Buffer
Volume:	2 vials (50 µL) negative lysate 2 vials (50 µL) positive lysate
Preparation:	Following cell treatment, NIH3T3 cell lysates were prepared on ice in MSD Complete Tris Lysis Buffer. Cell debris was cleared by centrifugation.
Storage:	Lysates should be stored at -80°C. Lysates will retain approximately 90% of activity after a single round of freeze thaw if handled properly (thawed on ice and immediately refrozen in smaller aliquots).
Quality Control:	Lysates have been tested for performance in Western Blot and MSD MULTI-ARRAY® Assays.

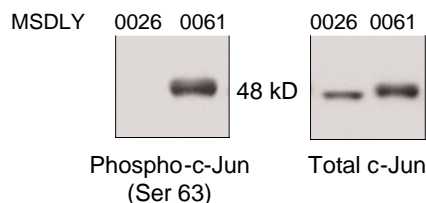
MSD MULTI-ARRAY Assay Results

The figure below illustrates a typical titration for MSDLY0026 (pc-Jun negative) and MSDLY0061 (pc-Jun positive) cell lysates with the MSD MULTI-ARRAY Phospho-c-Jun (Ser 63) Whole Cell Lysate Kit. The results are presented as a ratio of the signals obtained with pc-Jun positive and pc-Jun negative lysates. The phospho-c-Jun signal ratios increase with the amount of lysate. The representative results shown below are for demonstration purposes only and individual results may vary depending upon experimental application.



Traditional Western Blot Results

MSDLY0026 and MSDLY0061 whole cell lysates (20 µg each) were analyzed by Western Blot with phospho specific c-Jun and total c-Jun antibodies.



FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

20201-v1-2010Feb